

# Thermo Scientific Orion pH Electrode Storage Sleeve

The Thermo Scientific Orion pH electrode storage sleeve, 810017, is intended for use with all pH electrodes that are 120 mm or less in length that have a diameter of 12 mm. The storage sleeve provides protection for your electrode when not in use preventing breakage of the glass and also conditions the pH bulb so the electrode is ready to use when removing it from the storage sleeve.

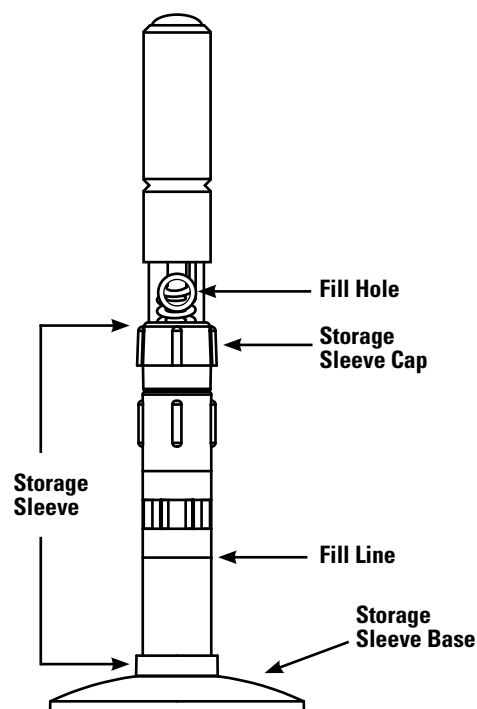
Follow the steps below to correctly install or remove your electrode from the storage sleeve.

## Installing the Electrode

1. Ensuring that the storage sleeve is in the upright position, add pH electrode storage solution in the sleeve up to the fill mark on the sleeve. It is recommended to use Orion 910001 for AgCl pH electrodes and 810001 for ROSS pH electrodes
2. Loosen the cap slightly to allow the electrode to gently slide in. The cap does not need to be removed to insert or remove the electrode. For low maintenance electrodes without a fill hole, the electrode should be fully placed in the sleeve so the cap touches the top of the storage sleeve. If it is a refillable electrode with a fill hole, the fill hole should be visible above the cap. Otherwise the fill hole should be sealed with tape to prevent storage solution from entering the electrode through the fill hole and the electrode can be placed fully inside the storage sleeve. Finger tighten the cap and place the storage sleeve in the base.

## Removing the Electrode

1. Slightly loosen the red storage sleeve cap.
2. With the storage sleeve in the upright position, gently remove the electrode from the sleeve realizing the pH bulb is delicate.
3. Store the sleeve and base upright to prevent spilling of the storage solution. If the electrode will be out of the storage sleeve for an extended period, it is best to remove the solution from the sleeve and replace with fresh storage solution when needed again.



Thermo Fisher Scientific  
Water Analysis Instruments  
Chelmsford, MA USA  
Quality Management System  
Registered to ISO 9001

[thermoscientific.com/water](http://thermoscientific.com/water)

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. & its subsidiaries

### Water Analysis Instruments

**North America**  
Toll Free: 1-800-225-1480  
Tel: 1-978-232-6000  
info.water@thermo.com

**Netherlands**  
Tel: (31) 033-2463887  
info.water.uk@thermo.com

**China**  
Tel: (86) 21-68654588  
wai.asia@thermofisher.com

**India**  
Tel: (91) 22-4157-8800  
wai.asia@thermofisher.com

**Singapore**  
Tel: (65) 6778-6876  
wai.asia@thermofisher.com

**Japan**  
Tel: (81) 045-453-9175  
wai.asia@thermofisher.com

**Australia**  
Tel: (613) 9757-4300  
in Australia (1300) 735-295  
InfoWaterAU@thermofisher.com

**Thermo**  
SCIENTIFIC

Part of Thermo Fisher Scientific